

AMPIVIEW[®] DNP (anti-rabbit) Detection AP-Red Kit

Complete staining kit with highest sensitivity, lowest background

AMPIVIEW[®] DNP (anti-rabbit) Detection AP-Red Kit is intended for AMPIVIEW[®] RNA probes ISH staining. The kit includes AMPIVIEW[®] Rabbit Anti-DNP linker, POLYVIEW[®] PLUS AP IHC reagent, HIGHDEF[®] chromogen and counterstain, and it is suitable visualization of the spatial localization of target genes in formalin-fixed paraffin-embedded (FFPE) tissue specimens without disrupting morphology.

Ordering Information

[Order Online »](#)

ENZ-ACC176-0020	20 tests
-----------------	----------

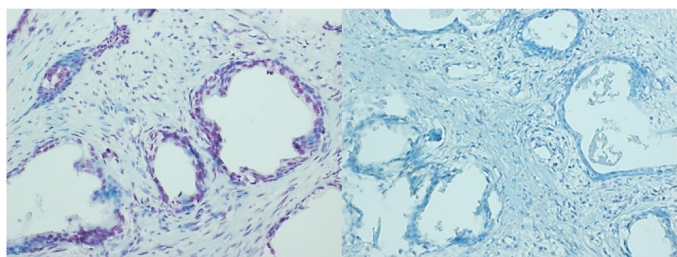
Manuals, SDS & CofA

[View Online »](#)

- Highest sensitivity, complete staining kit for DNP-labeled AMPIVIEW[®] RNA Probes
- POLYVIEW[®] PLUS reagents' biotin free nanopolymer detection circumvents endogenous background while providing specific antigen localization
- HIGHDEF[®] chromogens provides high intensity color development to deliver sharp, crisp staining
- Compatible with both manual and automated platforms

A.

B.



PTEN detected in prostate cancer tissue with A. AMPIVIEW® PTEN (AS) DNP RNA Probes (red) and B. AMPIVIEW® NSP DNP RNA Probes (no signal). Tissue samples were pre-treated with antigen retrieval reagent, pH 9 (ENZ-ACC113), Proteinase K (ENZ-33801), and AMPIVIEW® PTEN (AS) DNP RNA probes (ENZ-GEN131) were amplified with AMPIVIEW® anti-DNP Linker (RTU) (ENZ-ACC928) and detected with POLYVIEW® PLUS (anti-rabbit) AP (ENZ-ACC110), combined with HIGHDEF® Red AP chromogen kit (ENZ-ACC158) and hematoxylin (ENZ-ACC106).

Handling & Storage

Handling Do not freeze.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application IHC, ISH (in situ hybridization)

Contents POLYVIEW[®] PLUS AP (Anti-Rabbit) Reagent (Ready-To-Use) (ENZ-ACC110), HIGHDEF[®] Red Chromogen/Substrate, HIGHDEF[®] Hematoxylin (ENZ-ACC106), Antigen Retrieval Reagent, pH 9 (10X) (ENZ-ACC113), AMPIVIEW[®] Rabbit anti-DNP linker (ENZ-ABS928), Proteinase K (ENZ-33801), AMPIVIEW[®] Wash Buffer 1, AMPIVIEW[®] Wash Buffer 2

Technical Info / Product Notes

AMPIVIEW[®] DNP (anti-rabbit) Detection AP-Red Kit is a convenient detection kit for AMPIVIEW[®] RNA probes with reagents from antigen retrieval to counterstain. AMPIVIEW[®] RNA Probes are uniquely designed with the precision of targeted, sequence specific RNA, powered by Enzo's LoopRNA[™] ISH technology to deliver superior sensitivity. Sample specimens are pre-treated with antigen retrieval and Proteinase K treatment before dinitrophenol (DNP)-labeled AMPIVIEW[®] RNA probes are amplified with AMPIVIEW[®] anti-DNP linker and detected with POLYVIEW[®] PLUS detection reagents combined with HIGHDEF[®] chromogens and counterstain. Spatial biology of target genes can be visualized within its morphological context under the light microscope.

Last modified: January 13, 2025



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)